

January 5, 2022

NutriFreez® D10 Cryopreservation Medium

Scientific References

- **Pluripotent Stem Cells**
- **Mesenchymal Stem Cells**
- **Peripheral Blood Mononuclear Cells (PBMC)**
- **Immune cells**
- **Other Cell Lines**

Pluripotent Stem Cells

1. L. E. Wiersma, et al. **Large-Scale Engineering and Cryopreservation of hiPSC-Derived Nephron Sheets.** *Stem Cell Research & Therapy*, 2021. DOI: <https://doi.org/10.21203/rs.3.rs-1116352/v1>
2. M. Breunig, et al. **Differentiation of human pluripotent stem cells into pancreatic duct-like organoids.** *STAR Protocols*, 2021, <https://doi.org/10.1016/j.xpro.2021.100913>.
3. S.E. Strawbridge, et al. **Deriving Human Naïve Embryonic Stem Cell Lines from Donated Supernumerary Embryos Using Physical Distancing and Signal Inhibition.** *Methods in Molecular Biology, Humana*, https://doi.org/10.1007/978-1-0716-1908-7_1
4. E. Balbasi, et al. **Mouse Embryonic Stem Cell Culture in Serum-Containing or 2i Conditions.** *Methods in Molecular Biology.* Springer, 2021. https://doi.org/10.1007/978-1-0716-1908-7_1
5. D. Khan, et al. **Testing the Stability of Drug Resistance on Cryopreserved, Gene-Engineered Human Induced Pluripotent Stem Cells.** *Pharmaceuticals* 2021, <https://doi.org/10.3390/ph14090919>
6. B. Liu, et al. **Chemically defined and xeno-free culture condition for human extended pluripotent stem cells.** *Nat Commun* (2021). <https://doi.org/10.1038/s41467-021-23320-8>
7. R.R. Annand, **Generation of Human iPSCs by Reprogramming with the Unmodified Synthetic mRNA.** *Nuclear Reprogramming.* Part of *Methods in Molecular Biology*, (2021) vol 2239. Humana, New York, NY. https://doi.org/10.1007/978-1-0716-1084-8_11
8. J. Selberg, et al. **Machine Learning-Driven Bioelectronics for Closed-Loop Control of Cells.** *Adv. Intell. Syst.* 2020, DOI: 10.1002/aisy.202000140
9. Z. Danhong, et al. **Method of Cryopreservation of Stem Cell-Derived Retinal Pigment Epithelial Cells on Polymeric Substrate.** Application Number: 16/658457, Publication Date: 02/13/2020
10. M. G. Hernandez, et al. **AAV-Mediated Gene Delivery to 3D Retinal Organoids Derived from Human Induced Pluripotent Stem Cells.** *Int. J. Mol. Sci.* 2020, 21, 994; doi:10.3390/ijms21030994
11. S. Kumar, et al. **Highly efficient induced pluripotent stem cell reprogramming of cryopreserved lymphoblastoid cell lines.** *Journal of Biological Methods*, 2020, Vol. 7(1), e124, DOI: 10.14440/jbm.2020.296
12. L. Kuznitsov-Yanovsky, et al. **Modeling FXS: Human Pluripotent Stem Cells and In Vitro Neural Differentiation.** In: Ben-Yosef D., Mayshar Y. (eds) *Fragile-X Syndrome.* *Methods in Molecular Biology*, vol 1942. Humana Press, New York, NY (2019) https://doi.org/10.1007/978-1-4939-9080-1_8
13. G. Gagliardi, et al. **Characterization and Transplantation of CD73-Positive Photoreceptors Isolated from Human iPSC-Derived Retinal Organoids.** *Stem Cell Reports j* Vol. 11 j 665-680 j September 11, 2018
14. M. Amit & J. Itskovitz-Eldor, **Single cells pluripotent stem cells in a suspension culture.** US Patent App. 16/055,110, 2018
15. G. Guo, et al. **Epigenetic resetting of human pluripotency.** *bioRxiv* (2017): 146712.
16. S. Reichman, et al. **Generation of Storable Retinal Organoids and Retinal Pigmented Epithelium from Adherent Human iPSCs in Xeno-Free and Feeder-Free Conditions.** *STEM CELLS* (2017): 1176-1188.
17. F. Pistollato, et al. **Protocol for the Differentiation of Human Induced Pluripotent Stem Cells into Mixed Cultures of Neurons and Glia for Neurotoxicity Testing.** *JoVE (Journal of Visualized Experiments)* 124 (2017): e55702-e55702.
18. L. Tian, et al. **Efficient and controlled generation of 2D and 3D bile duct tissue from human pluripotent stem cell-derived spheroids.** *Stem Cell Reviews and Reports* 12.4 (2016): 500-508.
19. J. Chal, et al. **Generation of human muscle fibers and satellite-like cells from human pluripotent stem cells in vitro.** *Nature protocols* 11.10 (2016): 1833-1850.
20. M. S. Y Hiew. **Generation and Microrna Expression Profiling of Colorectal Cancer Cell-Derived Induced Pluripotent Cancer Cells.** Diss. UTAR, 2016.
21. N. Nishishita, et al. **An effective freezing/thawing method for human pluripotent stem cells cultured in chemically-defined and feeder-free conditions.** *American journal of stem cells* 4.1 (2015): 38.
22. Z. Danhong, et al. **Method of cryopreservation of stem cell-derived retinal pigment epithelial cells on polymeric substrate.** U.S. Patent Application No. 14/114,182, 2014
23. C. Martin, et al. **Highly Efficient microRNA-Enhanced mRNA Reprogramming of Diseased Human Fibroblasts in a Feeder-Free Culture System.** *STEMGENT Application Note*, 2015
24. C. Buensuceso, et al. **Induced pluripotent stem cells prepared from human kidney-derived cells.** US Patent 20140073049 A1, 2014
25. R. Sivapatham, X.Zeng. **Generation and Characterization of Patient-Specific Induced Pluripotent Stem Cell for Disease Modeling.** *Methods in Molecular Biology*, 2014
26. L. Tian, et al. **In Vitro Modeling of Alcohol-Induced Liver Injury Using Human-Induced Pluripotent Stem Cells.** *Methods in Molecular Biology*, 2014

Mesenchymal Stem Cells

1. N. Aydođdu, et al. **Isolation, Culture, Cryopreservation, and Preparation of Umbilical Cord-Derived Mesenchymal Stem Cells as a Final Cellular Product Under Good Manufacturing Practices-Compliant Conditions.** *Methods in Molecular Biology.* (2020) Springer, New York, NY. https://doi.org/10.1007/7651_2020_332
2. N. Aydođdu, et al. **Isolation, Culture, Cryopreservation, and Preparation of Skin-Derived Fibroblasts as a Final Cellular Product Under Good Manufacturing Practice-Compliant Conditions.** *Methods in Molecular Biology.* (2020) Springer, New York, NY. https://doi.org/10.1007/7651_2020_333
3. M. Paz Quesada, et al. **Safety and Biodistribution of Human Bone Marrow-Derived Mesenchymal Stromal Cells Injected Intrathecally in Non-Obese Diabetic Severe Combined Immunodeficiency Mice: Preclinical Study.** *Tissue Eng Regen Med*, 03 July 2019, <https://doi.org/10.1007/s13770-019-00202-1>
4. A.M. Lyness, et al. **Evaluation of a Vial Adaptor to Ensure Safe and Efficient Needle-Free Transfer of Cells Post-Cryopreservation in Preparation for Drug Delivery.** *Cytotherapy*, Volume 21, Issue 5, Supplement, May 2019, Page S38
5. M. Salkhordeh, et al. **Evaluation of Different Cryopreservation Agents for Mesenchymal Stem Cell as Final Study Product.** *Cytotherapy*, Volume 20, Issue 5, Supplement, May 2018, Page S50
6. G. Kundrotas, et al. **Identity, proliferation capacity, genomic stability and novel senescence markers of mesenchymal stem cells isolated from low volume of human bone marrow.** *Oncotarget* 7.10 (2016): 10788.
7. G. Kundrotas, **Investigation of human mesenchymal stem cell genetic characteristics to ensure therapy safety.** Diss. Vilnius University, 2016.

Peripheral Blood Mononuclear Cells (PBMC)

1. N. Milevoj, et. al. **A combination of electrochemotherapy, gene electrotransfer of plasmid encoding canine IL-12 and cyto-reductive surgery in the treatment of canine oral malignant melanoma.** *Research in Veterinary Science*, volume 122, February 2019, Pages 40-49
2. R. E. Burnham, et al. **Human serum albumin and chromatin condensation rescue ex vivo expanded $\gamma\delta$ T cells from the effects of cryopreservation.** *Cryobiology*, 2021, ISSN 0011-2240, <https://doi.org/10.1016/j.cryobiol.2021.01.011>.

Immune cells

1. L. Xiao, et al. **Spectral Markers for T Cell Death and Apoptosis—A Pilot Study on Cell Therapy Drug Product Characterization Using Raman Spectroscopy.** *Journal of Pharmaceutical Sciences*, 2021, <https://doi.org/10.1016/j.xphs.2021.08.005>
2. R. E. Burnham, et al. **Human serum albumin and chromatin condensation rescue ex vivo expanded $\gamma\delta$ T cells from the effects of cryopreservation.** *Cryobiology*, 2021, ISSN 0011-2240, <https://doi.org/10.1016/j.cryobiol.2021.01.011>.
3. N. Milevoj et. al. **A combination of electrochemotherapy, gene electrotransfer of plasmid encoding canine IL-12 and cyto-reductive surgery in the treatment of canine oral malignant melanoma.** *Research in Veterinary Science*, volume 122, February 2019, Pages 40-49

Other Cell Lines

1. D. Pamies, et al. **Guidance document on Good Cell and Tissue Culture Practice 2.0 (GCCP 2.0)**. ALTEX. 2021 Dec 9. doi: 10.14573/altex.2111011.
2. Y. Jing-Yan, et al. **Retinal Protection by Sustained Nanoparticle Delivery of Oncostatin M and Ciliary Neurotrophic Factor Into Rodent Models of Retinal Degeneration**. Translational Vision Science & Technology August 2021, Vol.10, 6. doi:https://doi.org/10.1167/tvst.10.9.6
3. S. Reichman, et al. **Compositions and Methods for Efficient Amplification of Retinal Progenitors Cells**. US Patent App. 16/976581, 01/07/2021
4. L. Klipfel, et al. **A Splice Variant in SLC16A8 Gene Leads to Lactate Transport Deficit in Human iPSC-Derived Retinal Pigment Epithelial Cells**. Cells 2021,10, 179. https://doi.org/10.3390/cells10010179
5. X. Jiang, et al. **Mitochondria Dynamically Transplant into Cells in Vitro and in Mice and Rescue Aerobic Respiration of Mitochondrial DNA-Depleted Motor Neuron NSC-34**. Journal of Biomedical Science and Engineering, 13, 203-221 (2020). doi: 10.4236/jbise.2020.139019.
6. A. C. Brown, et al. **Long-Term Culture of Nephron Progenitor Cells Ex Vivo**. In: Vainio S. (eds) Kidney Organogenesis. Methods in Molecular Biology, vol 1926. Humana Press, New York, NY
7. E. Nigro, et al. **An Antitumor Cellular Vaccine Based on a Mini-Membrane IgE**. J Immunol 188:103-110 (2012).
8. G. Di Lullo, et al. **The combination of marker gene swapping and fluorescence-activated cell sorting improves the efficiency of recombinant modified vaccinia virus Ankara vaccine production for human use**. J. of Virological Methods, 163, Issue 2: 195-204 (2010)
9. M. Amit, & J. Itskovitz-Eldor. **Novel methods and culture media for culturing pluripotent stem cells**. U.S. Patent Application 13/821,244.
10. E. De Falco, et al. **A standardized laboratory and surgical method for in vitro culture isolation and expansion of primary human Tenon's fibroblasts**. Cell and tissue banking 14.2 (2013): 277-287.

Germany

Sartorius Stedim Biotech
GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North
America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178